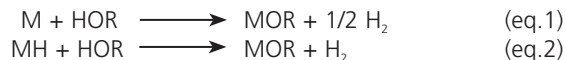


**AL-123**

Cat. No. Z516198

Technical Bulletin**Quantitative Analysis of Active Metals and Metal Hydrides via Gas Evolution**

Active metals (M) and metal hydrides (MH) (including boron hydrides) react rapidly and quantitatively with suitable hydrolytic solvents to produce hydrogen gas (eqs. 1 and 2).



The direct measurement of the volume of hydrogen gas produced provides a convenient and highly accurate method for the determination of concentration or purity of these reagents. Metals and metal hydrides that can be analyzed in this manner include sodium and lithium as dispersions in mineral oil, boranetetrahydrofuran and borane-methyl sulfide complexes, 9 BBN, K-Selectride®, lithium aluminum hydride, diisobutylaluminum hydride. Several reviews regarding the use and handling of boron hydrides and other active metals and metal hydrides are available.¹⁻⁷

A simple Aldrich Kit for measuring the volume of hydrogen gas is illustrated in Fig.1. This apparatus can be assembled from readily available laboratory equipment.

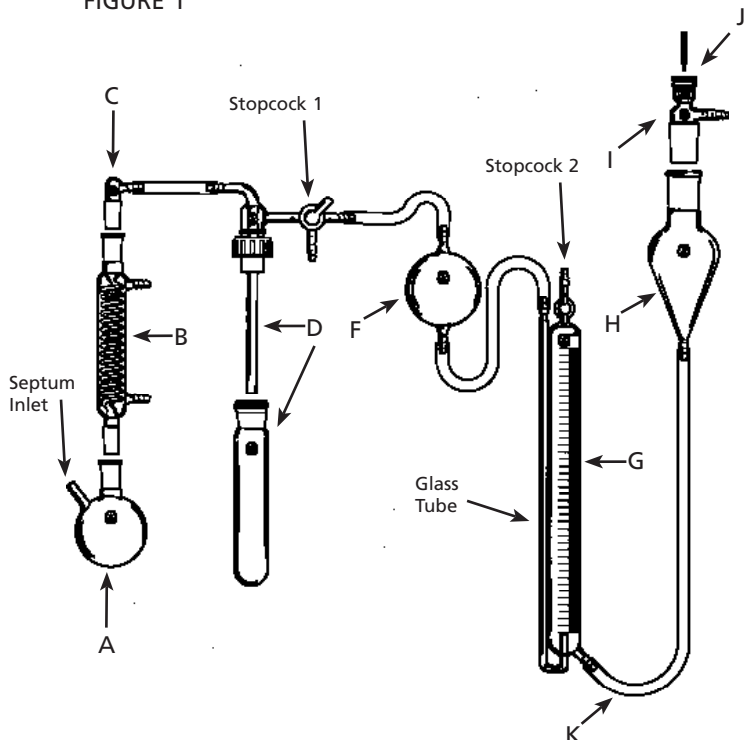
PROCEDURES**A. DETERMINATION OF HYDRIDE CONCENTRATION OF A SOLUTION**

The following procedure for determining the hydride concentration of a $BH_3 \cdot THF$ solution is representative. The gas burette is assembled as illustrated in Fig. 1 and the septum-inlet flask A is charged with 50ml of glycerol and 50ml of water. Cooling water is run through the Graham condenser B and trap D is cooled in a dry-ice/acetone bath. With stopcocks 1 and 2 both open to the atmosphere, the leveling bulb H is adjusted to give a zero reading for the level of distilled water in the gas burette. Stopcock 2 is then closed and stopcock 1 is turned so that the septum-inlet flask and the graduated burette form a closed system. An accurately measured aliquot of the $BH_3 \cdot THF$ solution (2.00ml of a 1M solution generates ca. 150ml of hydrogen) is added slowly through the septum-inlet to the stirred glycerol/water solution using a hypodermic syringe. 4,5 Hydrolysis of the $BH_3 \cdot THF$ solution is extremely rapid, being complete in a few seconds.

B. HYDRIDE DETERMINATION OF SOLID INORGANIC HYDRIDES OR HYDROGEN-GENERATING METALS

Assemble the kit as in Fig. 1, without the septum-inlet flask. Turn on the cooling water and place cold trap D in a dry ice/acetone bath. With stopcocks 1 and 2 open to the atmosphere, adjust the leveling bulb H to give a zero reading.

FIGURE 1



- A. Z101265 Septum-inlet flask, 250mL
- B. Z531413 Graham condenser, 300mm
- C. Z137057 Vacuum adapter
- D. Z551724 Cold-trap top with stopcock
- Z551732 Cold-trap bottom (dry ice/acetone)
- F. Z551740 Trap flask, 250mL
- G. Z551759 Graduated burette, 250mL with 1mL graduations
- H. Z551767 Leveling bulb, 250mL
- I. Z531782 Thermometer adapter
- J. Z253065 Non-mercury thermometer, total immersion, 0 to 230°
- F
- K. Z280380 PVC flexible tubing

Inside a glove bag and under nitrogen or argon, transfer 3-6 meq of sample to a septum-topped, flushed, tared 100ml septum-inlet-equipped flask containing a stirring bar. (An antistatic alpha-emitter plate held next to the neck of the flask during transfer controls loss of sample onto the neck of the flask.) Reseal the flask and reweigh. Working quickly, attach the sample flask in position under the condenser. Add 10ml of anhydrous solvent (which has been stored over calcium hydride) through the **septum inlet** to slurry the solid. Close stopcock **2**; turn stopcock **1** so that the septum inlet flask and graduated buret form a closed system. Slowly inject 2.0ml of hydrolysis solution. In most cases, hydrolysis is complete in less than 2 minutes.

When hydrolysis is complete, the water level in **glass tube** will remain constant at a level considerably below the water level in the buret. To obtain an accurate reading of the volume of hydrogen gas produced, the water level in the **glass tube** must be coplanar with the water level in the graduated buret **G**. Consequently, more hydrogen gas must be forced over into the burette to raise the water level in the side tube until it corresponds to the water level in the buret. This is easily accomplished by quickly lowering the leveling bulb **H** to a point below the outlet of the side tube in the buret and then raising the leveling bulb to its original position. The lowering and raising of the leveling bulb are repeated until the water levels in the side tube, burette and leveling bulb are all coplanar. The final burette reading is then taken as illustrated in **Fig. 1**.

The analysis is repeated until three separate analyses vary by no more than 1ml. Occasionally, the initial analysis will be highly inaccurate and this result should be discarded. When consistent measurements are obtained for the gas volume, the molarity of the borane solution is calculated using the following equation:

$$\text{Borane molarity} = \frac{(P_1 - P_2) (273) (V_1 - V_2)}{(760) (T) (22.4) (V_2) (n)}$$

Where:

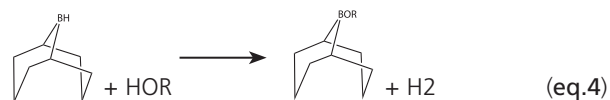
- P_1 = observed pressure (mm Hg)
- P_2 = vapor pressure (mm Hg) of water at T (see **Table 1**)
- V_1 = volume (ml) of hydrogen gas evolved
- V_2 = volume (ml) of borane solution injected into hydrolyzing solution
- T = observed temperature ($^{\circ}\text{K}$)
- n = number of moles of hydrogen produced per mole of sample; The value of n is 3, for example, for $\text{BH}_3 \cdot \text{THF}$

since

3 moles of hydrogen are generated from one mole of $\text{BH}_3 \cdot \text{THF}$ solution at standard temperature and pressure based on the stoichiometry shown in **eq. 3**.



For 9-BBN, K-Selectride[®] and Super-Hydride[®], the value of n is 1, based on the stoichiometry shown in **eqs. 4 and 5**.



The procedure described for analysis of the BH_3/THF solution can be easily adapted to the analysis of solutions of other active hydrides by changing the hydrolysis solution and conditions. **Table 2** is a summary of the conditions necessary for the hydrolysis of various active hydride solutions.

The activity or purity of a solid compound is calculated using the following equation:

$$\% \text{MH}_n \text{ (or \%M)} = \frac{(P_1 - P_2) (273) (V_1 - V_3) (\text{MW}) (100)}{(760) (T) (22.4) (x) (n)}$$

Where:

- V_3 = volume of hydrolyzing solution added
- MW = molecular weight
- x = weight of sample in mg

Table 3 is a summary of solvents and hydrolyzing solutions necessary for the hydrolysis of various solid metal hydrides and hydrogen-generating metals.

C. DETERMINATION OF PURITY OF SOLID OR LIQUID BORON HYDRIDES

The simplest procedure for solid or liquid boron hydrides involves first dissolving an accurately weighed quantity of pure boron hydride in an appropriate dry solvent (e.g., THF) in a volumetric flask. Dilution to a known volume gives a standard solution. The theoretical molarity can then be calculated and compared with the experimental molarity determined by hydrogen evolution upon hydrolysis. **Table 4** summarizes the conditions necessary for the hydrolysis of various solid or liquid boranes dissolved in THF.

Borane-*tert*-butylamine, borane-dimethylamine, boranemorpholine, borane-pyridine, borane-triethylamine, and borane-trimethylamine complexes are all essentially unreactive toward hydrolytic solvents at neutral pH. These amine-boranes are hydrolyzed by aqueous mineral acids (**eq. 6**). However, the reaction is generally too slow to be used conveniently for quantitative analysis via hydrogen evolution; e.g., a period of 6 hours is necessary for the complete hydrolysis of borane-triethylamine by 1M HCl in water/ethylene glycol solution.⁵ These hydrolytically stable borane complexes are more conveniently analyzed by IR, NMR, GC or elemental analysis.

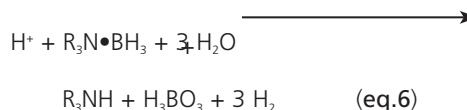


TABLE 1
VAPOR PRESSURE (MM HG) OF WATER*

Temp. °C	0.0	0.2	0.4	0.6	0.8
19	16.477	16.685	16.894	17.105	17.319
20	17.535	17.753	17.974	18.197	18.422
21	18.650	18.880	19.113	19.349	19.587
22	19.827	20.070	20.316	20.565	20.815
23	21.068	21.324	21.583	21.845	22.110
24	22.377	22.648	22.922	23.198	23.476
25	23.756	24.039	24.326	24.617	24.912

*Reprinted with permission from *Handbook of Chemistry and Physics*, 67th ed.; Weast, S.C., Ed; CRC Press, Inc: Boca Raton, 1986; p D-143. Copyright by CRC Press, Inc.

TABLE 2
CONDITIONS FOR THE HYDROLYSIS OF ACTIVE METAL-HYDRIDE SOLUTIONS

Catalog No.	Description	Solution ^a	Hydrolysis Rate ^b
229024	R-Alpine-Hydride [®]	G-W-THF	R
237728	S-Alpine-Hydride [®]	G-W-THF	R
193852	9-BBN, 0.5M in hexanes	M-THF	S
151076	9-BBN, 0.5M in THF	M-THF	S
258830	9-BBN-pyridine, 0.5M in diethyl ether	1	R
192112	Borane-methyl sulfide complex, 2.0M in diethyl ether	G-W-M	R
192120	Borane-methyl sulfide complex, 2.0M in THF	G-W-M	R
194824	Borane-methyl sulfide complex, 2.0M in toluene	G-W-M	R
176192	Borane-tetrahydrofuran complex, 1.0M in THF	G-W	R
225762	Catecholborane, 1.0M in THF	G-W-THF	R
214949	Diisobutylaluminum hydride, 1.0M in cyclohexane	2	R
190306	Diisobutylaluminum hydride, 1.0M in hexanes	2	R
214973	Diisobutylaluminum hydride, 1.0M in dichloromethane	2	R
214981	Diisobutylaluminum hydride, 1.0M in THF	2	R
215007	Diisobutylaluminum hydride, 1.0M in toluene	2	R
220760	K-Selectride [®]	G-W-THF	R
220779	KS-Selectride [®]	G-W-THF	R
212792	Lithium aluminum hydride, 1.0M in diethyl ether	3	R
236055	Lithium aluminum hydride, 0.5M in diglyme	3	R
212776	Lithium aluminum hydride, 1.0M in THF	3	R
243949	Lithium aluminum hydride bis(tetrahydrofuran), 1.0M in toluene	3	R
230200	Lithium borohydride, 2.0M in THF	3	R
244538	Lithium thexylimonylborohydride, 0.5M in THF	G-W-THF	R
178497	L-Selectride [®]	G-W-THF	R
225924	LS-Selectride [®]	G-W-THF	R
213403	N-Selectride [®]	G-W-THF	R
213438	Potassium triethylborohydride, 1.0M in THF	G-W-THF	R
197890	Potassium triisopropoxyborohydride, 1.0M in THF	G-W-THF	R
243922	1-Pyrrolylborane, 1.0M in THF	G-W-THF	R
200972	Sodium borohydride, 0.5M in diglyme	HCl-EG-W	R
215511	Sodium borohydride, 2.0M in triglyme	HCl-EG-W	R
227315	Sodium triethylborohydride, 1.0M in diethyl ether	G-W-THF	R
213411	Sodium triethylborohydride, 1.0M in THF	G-W-THF	R
227307	Sodium triethylborohydride, 1.0M in toluene	G-W-THF	R
180866	Super-Deuteride [®]	G-W-THF	R
179728	Super-Hydride [®]	G-W-THF	R

^a Hydrolysis solutions prepared by combining equal volumes of the solvents given where G = glycerol, W = water, M = methanol, EG = ethylene glycol, THF = tetrahydrofuran and HCl = 6M hydrochloric acid.

^b R = Rapid, hydrolysis complete in < 5 minutes at 20-25°. S = slow, hydrolysis requires 15-20 minutes for completion at 25°.

1 = Conc. HCl and methanol 1:1 cooled in ice bath

2 = Conc. HCl, water, hexanes 1:1:4 cooled in ice bath

3 = 2N HCl cooled in ice bath

HANDLING INSTRUCTIONS

The reagents discussed in this bulletin are air- and /or moisture sensitive. Most of the products are supplied in our Sure/Seal™, Sure/Pac™ or Kilo-Lab™ containers for extra protection. The products are to be handled, and the procedures outlined here performed only by technically qualified persons. For information on the techniques and equipment for handling these products safely, a number of Technical Information Bulletins⁴⁻⁸ is available from Aldrich. The bulletins are available free upon request.

REFERENCES AND NOTES

- 1 Brown, H.C.; Kramer, G.W.; Levy, A.B.; Midland, M.M. *Organic Synthesis via Boranes*; John Wiley and Sons, Inc.: New York, 1973; pp 21-244 (Aldrich Catalog No. Z101443).
- 2 Murray, L.I. "A Study of the Amine-boranes", Ph.D. Thesis, Purdue University, Lafayette, 1963.

- 3 *Handling Air-Sensitive Reagents*, Aldrich Technical Information Bulletin No. AL-134.
- 4 *Equipment for Handling Air-Sensitive Reagents*, Aldrich Technical Information Bulletin No. AL-135.
- 5 *The Aldrich Sure/Pac™ Cylinder Packaging System and Recommended Transfer Procedures*, Aldrich Technical Information Bulletin No. AL-136.
- 6 *Aldrich Kilo-Lab™ Cylinder Packaging System and Recommended Transfer Procedures*, Aldrich Technical Information Bulletin No. AL 149.
- 7 *Equipment for Transfer of Liquids from Aldrich Kilo-Lab™ Cylinders*, Aldrich Technical Information Bulletin No. AL-150.
- 8 Normally, a small amount of cupric nitrate is dissolved in the water to prevent algae growth.

™ Trademark of Aldrich Chemical Company, Inc.

® Registered trademark of Aldrich Chemical Company, Inc.

**TABLE 3
CONDITIONS FOR HYDROLYSIS OF SOLID METAL HYDRIDES AND ACTIVE METALS**

Catalog No.	Description	Solution ^a	Solution ^b
208027	Calcium hydride, -4+40 mesh	T	W
213268	Calcium hydride, -40 mesh	T	W
213322	Calcium hydride, -1+4 mesh	T	W
248819	Lithium, 25 wt % in mineral oil, high Na	THF	I-M
248827	Lithium, 25 wt % in mineral oil, low Na	THF	I-M
199877	Lithium aluminum hydride	T	W
201049	Lithium hydride	T	W
215813	Potassium hydride, 35% in mineral oil	THF	THF-W
217123	Sodium, 40 wt % in light oil	THF	THF-W
217115	Sodium, 40 wt % in odorless mineral spirits	THF	THF-W
198072	Sodium borohydride	THF	HCl-EG-W
223441	Sodium hydride	THF	I-M
199230	Sodium hydride, 60 wt % in mineral oil	T	W
253995	Sodium hydride, 80 wt % in mineral oil	T	W

^a Solvent is 10ml of either toluene (T) or tetrahydrofuran (THF)

^b Hydrolysis solution mixtures are prepared by combining equal volumes of solvents where W = water, I = isopropanol, M = methanol, HCl = 6N hydrochloric acid and EG = ethylene glycol

**TABLE 4
CONDITIONS FOR THE HYDROLYSIS OF PURE BORON HYDRIDES**

Catalog No.	Description	Solution ^a	Rate ^b
178713	9-BBN	M-THF	S
179043	Borane-N,N-diethylaniline	G-W-THF	S
253111	Borane-N,N-diisopropylethylamine	EG-HCl	R
179825	Borane-methyl sulfide	G-W-M	R
224901	Borane-1,4-oxathiane	G-W-M	R
179833	Borane-4-phenylmorpholine	G-W	R
179752	Borane-pyridine	H ₂ SO ₄ -W-EG	R

^a Hydrolysis solutions prepared by combining equal volumes of the solvents given; where G = glycerol, W = water, M = methanol, EG = ethylene glycol, THF = tetrahydrofuran and HCl = 6N hydrochloric acid

^b R = Rapid, hydrolysis complete in < 5 minutes at 20-25°. S = Slow, hydrolysis requires 15-20minutes for completion at 25°.

PRECISION SEAL RUBBER SEPTA

Engineered for a precision fit in standard taper glassware joints and tubes. Manufactured under "White Room" conditions, from one certified raw material formulation for absolute consistency in all sizes, from lot to lot.

- Longer, tapered plug design
- Increases the sealing surface 30% while improving glass to rubber contact up to 80%
- Smooth, radial bottom edge
- Flexible sleeve

Description	<i>White</i>	<i>Red</i>
	Cat. No.	Cat. No
Fits 7mm O.D. tube	Z553905-10EA Z553905-100EA	Z554022-10EA Z554022-100EA
Fits 8mm O.D. tube	Z553913-10EA Z553913-100EA	Z554030-10EA Z554030-100EA
Fits 10mm O.D. tube	Z553921-10EA Z553921-100EA	Z554049-10EA Z554049-100EA
Fits 13mm O.D. tube	Z553948-10EA Z553948-100EA	Z554057-10EA Z554057-100EA
Fits 10/30 joint	Z553956-10EA Z553956-100EA	Z554065-10EA Z554065-100EA
Fits 14/20 joint	Z553964-10EA Z553964-100EA	Z554073-10EA Z554073-100EA
Fits 19/22 joint	Z553972-10EA Z553972-100EA	Z554081-10EA Z554081-100EA
Fits 24/40 joint	Z553980-10EA Z553980-100EA	Z554103-10EA Z554103-100EA
Fits 29/42 joint	Z553999-10EA Z553999-100EA	Z554111-10EA Z554111-100EA



We are committed to the success of our Customers, Employees and Shareholders through leadership in Life Science, High Technology and Service.

© 2006 Sigma-Aldrich Co. Printed in the USA. Aldrich is a member of the Sigma-Aldrich family. Aldrich warrants that its products conform to the information contained in this and other Aldrich publications. Purchaser must determine the suitability of the product for its particular use. See reverse of invoice of packing slip for additional terms and conditions of sale.